

## Blocking Immobilized Label-free Plates

### Introduction

The following short (optional) protocol shows how to block EnSpire-LFB plates (label free enabled plates for biochemical assays on the EnSpire<sup>®</sup> Multimode Plate Reader) following protein immobilization protocols. This step is not required for overnight immobilizations, as the surface chemistry of the plate should be completely deactivated (hydrolyzed) following an overnight incubation.

### Materials

- Ethanolamine (Sigma-Aldrich<sup>®</sup> PN 398136)
- 1X Phosphate Buffered Saline (PBS) (Sigma-Aldrich<sup>®</sup> PN P-3813), or appropriate assay buffer

### Blocking Plates

To deactivate residual active groups on the plate's surface, briefly wash in a buffer containing amine. To do this:

- Prepare a solution containing 50-100 mM ethanolamine in PBS
- Remove the current solution from all wells
- Add 40  $\mu$ L ethanolamine containing 1X PBS to the wells
- Incubate for 5 min at room temperature
- Wash the plate again 4 times with 25  $\mu$ L PBS
- Leave 15  $\mu$ L PBS in each well

PerkinElmer, Inc.  
940 Winter Street  
Waltham, MA 02451 USA  
P: (800) 762-4000 or  
(+1) 203-925-4602  
[www.perkinelmer.com](http://www.perkinelmer.com)



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright ©2014, PerkinElmer, Inc. All rights reserved. PerkinElmer<sup>®</sup> is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

May 22 2014